

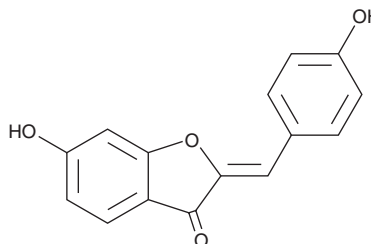
PRODUCT INFORMATION



Hispidol

Item No. 37373

CAS Registry No.: 5786-54-9
Formal Name: 6-hydroxy-2Z-[(4-hydroxyphenyl)methylene]-3(2H)-benzofuranone
Synonyms: 4',6-Dihydroxyaurone, (Z)-Hispidol
MF: C₁₅H₁₀O₄
FW: 254.2
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 380 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Hispidol is supplied as a solid. A stock solution may be made by dissolving the hispidol in the solvent of choice, which should be purged with an inert gas. Hispidol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hispidol in these solvents is approximately 10, 25, and 50 mg/ml, respectively.

Description

Hispidol is an aurone originally isolated from soybeans that has diverse biological activities.¹⁻⁴ It is an inhibitor of tyrosinase, monoamine oxidase A (MAO-A), and MAO-B (IC₅₀s= 38.4, 0.26, and 2.45 μM, respectively).^{1,3} Hispidol (50 μM) increases catalase and superoxide dismutase (SOD) activity in cell-free assays, as well as decreases intracellular reactive oxygen species (ROS) in, and increases the lifespan of, *C. elegans*.⁴ It decreases immobility time in forced swim and tail suspension tests in mice when administered at a dose of 2, 5, and 15 mg/kg.³

References

1. Okombi, S., Rival, D., Bonnet, S., *et al.* Discovery of benzylidenebenzofuran-3(2H)-one (aurones) as inhibitors of tyrosinase derived from human melanocytes. *J. Med. Chem.* **49**(1), 329-333 (2006).
2. Wong, E. Occurrence and biosynthesis of 4',6-dihydroxyaurone in soybean. *Phytochem.* **5**(3), 463-467 (1966).
3. Oh, J.M., Lee, H.-S., Baek, S.C., *et al.* Antidepressant-like activities of hispidol and decursin in mice and analysis of neurotransmitter monoamines. *Neurochem. Res.* **45**(8), 1930-1940 (2020).
4. Lim, H.J., Han, Y.T., Ahn, J.-H., *et al.* Longevity effects of hispidol in *Caenorhabditis elegans*. *Biofactors* **46**(6), 1041-1048 (2020).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/27/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM